

Data Sheet

TAF1 (1400-1518)

Human, recombinant, N-terminal HIS-tag

Catalog #: 31123

Lot #: 120827-D Conc.: 1.9 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Sako W, *et al.*, *Neuroscience*. 2011 Aug **25**;189:100-7.
2. Zaborowska J, *et al.*, *Transcription*. 2012 Mar 1;**3(2)**: 92-104.

Description:

Human Bromodomain TAF1 (TBP-associated factor), amino-acids 1400 – 1518 (GenBank Accession No. NM_004606) with N-terminal HIS-tag, MW = 14.9 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

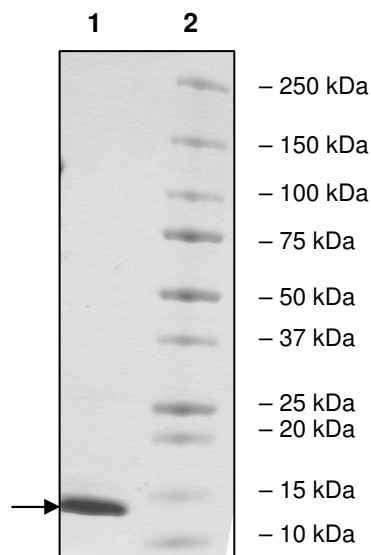
1.9 µg TAF1

Lane 2:

Protein Marker
Biorad (#161-0373)

MW: 14.9 kDa

Purity: ≥95%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Amino Acid Sequence:

MHHHHHHTDPMVTLSSILESIINDMRDLPNTYPFHTPVNAKVVVDYKIIITRPMDLQTLR
ENVRKRLYPSREEFHLELIVKNSATYNGPKHSLTQISQSMLDLCDEKLKEKEDKLA
RLEKAINP

Green: His-tag

Black: TAF1 (aa 1400-1518)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com